

VeriKine-HS Human Interferon Alpha All Subtype TCM ELISA Kit

PBL's VeriKine-HS Human Interferon Alpha All Subtype TCM ELISA Kit is specifically formulated to detect all human IFN-Alpha subtypes. This kit quantitates all human IFN alpha subtypes in tissue culture media (TCM) using a sandwich immunoassay format. The kit is based on an ELISA with biotinylated-detection antibody and streptavidin-conjugated horseradish peroxidase (HRP). Tetramethyl-benzidine (TMB) is the substrate. The assay is based on the international reference standard for human interferon alpha provided by the National Institutes of Health.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

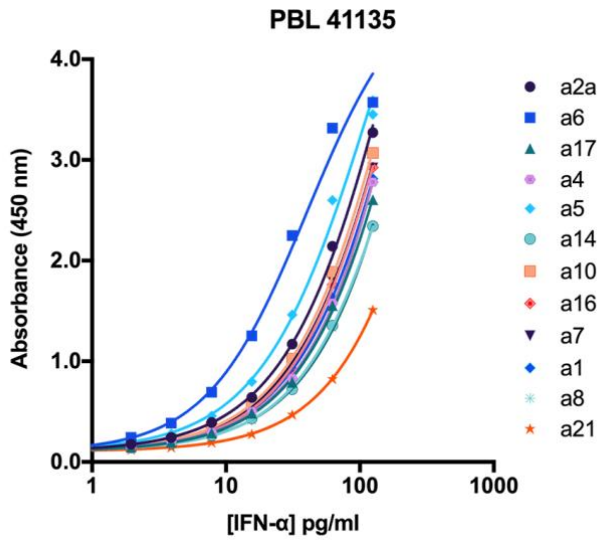
PBL-41135-1

96 wells

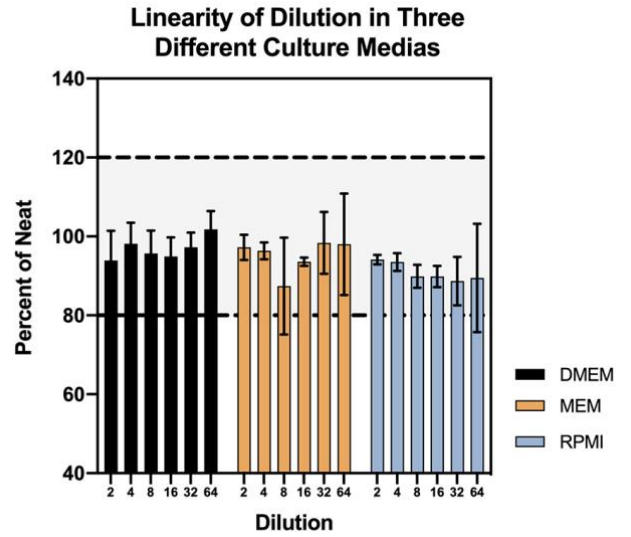
Manuals, SDS & CofA

[View Online »](#)

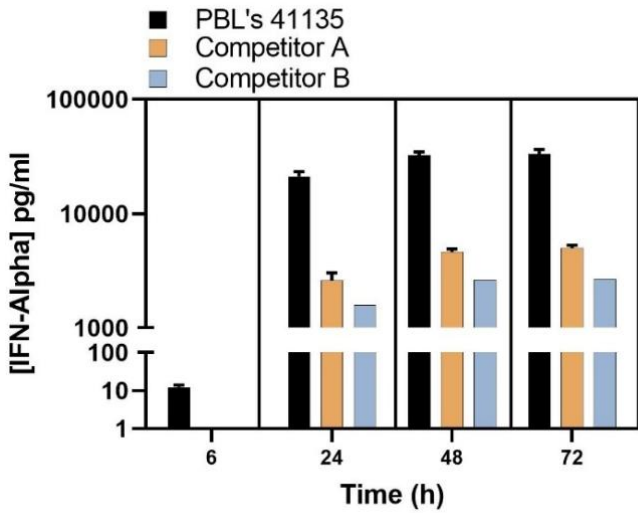
- Highly sensitive ELISA for global IFN alpha detection in tissue culture media (TCM)
- Low pg/ml detection of ALL human IFN- α subtypes
- Reproducible results with < 10% inter- and < 8% intra-assay CVs



Typical Human IFN-Alpha Subtype standard curves



Comparison of Endogenous Levels of IFN-Alpha quantified in Sendai Virus infected U937 Cells



Linearity of Dilution of Three Different Culture Medias

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN- α , Interferon- α

Application ELISA

Application Notes Quantitative determination of Hu-IFN alpha A, 2, A/D, B2, C, D, G, H, I, J1, K, 1, 4b and WA in tissue culture media

Assay Time 20 hours 30 minutes – 24 hours 30 minutes

Contents Pre-coated microtiter plate, Plate Sealers, Wash Solution Concentrate, Human Interferon Alpha Standard, 10,000 pg/ml, Assay Buffer, Standard Diluent, Antibody Diluent, Antibody Concentrate, HRP Conjugate Concentrate, HRP Diluent, TMB Substrate Solution, Stop Solution

Crossreactivity Cross-reacts with Cynomolgus/Rhesus IFN- α (13%). No cross-reactivity against human IFN- β , IFN- γ , or IFN- ω ; mouse or rat IFN- α , IFN- β or IFN- γ ; pig IFN- α ; or bovine IFN- τ .

Sensitivity 1.95 – 125 pg/ml

Species Reactivity Human



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)